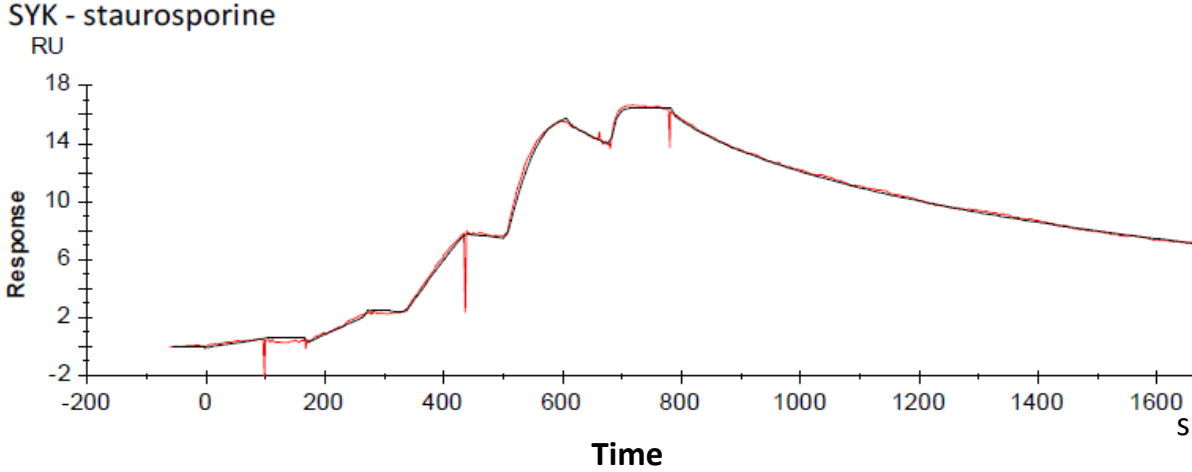


Binding kinetics of staurosporine on SYK



The data was generated and provided by
Netherlands Translational Research Center (NTRC)